

# PRODUCT INFORMATION

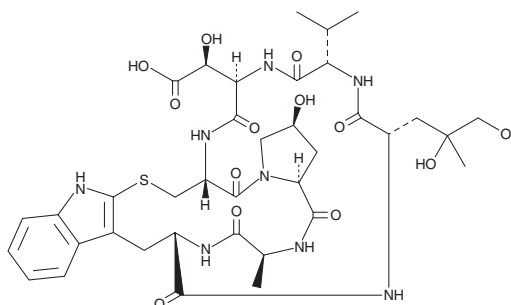


## Phalloidin

Item No. 18146

**CAS Registry No.:** 26645-35-2  
**Formal Name:** cyclic (2→6)-thioether-cyclo[L-alanyl-2-mercapto-L-tryptophyl-4,5-dihydroxy-L-leucyl-L-valyl-(3S)-3-hydroxy-D-α-aspartyl-L-cysteinyl-(4S)-4-hydroxy-L-prolyl]

**MF:** C<sub>37</sub>H<sub>50</sub>N<sub>8</sub>O<sub>13</sub>S  
**FW:** 846.9  
**Purity:** ≥90%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Phalloidin is supplied as a crystalline solid. A stock solution may be made by dissolving the phalloidin in the solvent of choice, which should be purged with an inert gas. Phalloidin is soluble in organic solvents such as methanol, DMSO, and dimethyl formamide. Phalloidin is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

### Description

Phalloidin is a natural mycotoxin first isolated from the death cap mushroom, *A. phalloides*. It binds filamentous actin (F-actin) like phalloidin (Item No. 18039), but contains a carboxyl group for coupling reactions.<sup>1,2</sup>

### References

1. Barak, L.S., Yocum, R.R., Nothnagel, E.A., *et al.* Fluorescence staining of the actin cytoskeleton in living cells with 7-nitrobenz-2-oxa-1,3-diazole-phalloidin. *Proc. Natl. Acad. Sci. USA* **77**(2), 980-984 (1980).
2. Barak, L.S., Yocum, R.R., and Webb, W.W. *In vivo* staining of cytoskeletal actin by autointernalization of nontoxic concentrations of nitrobenzoxadiazole-phalloidin. *J. Cell Biol.* **89**, 368-372 (1981).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM